

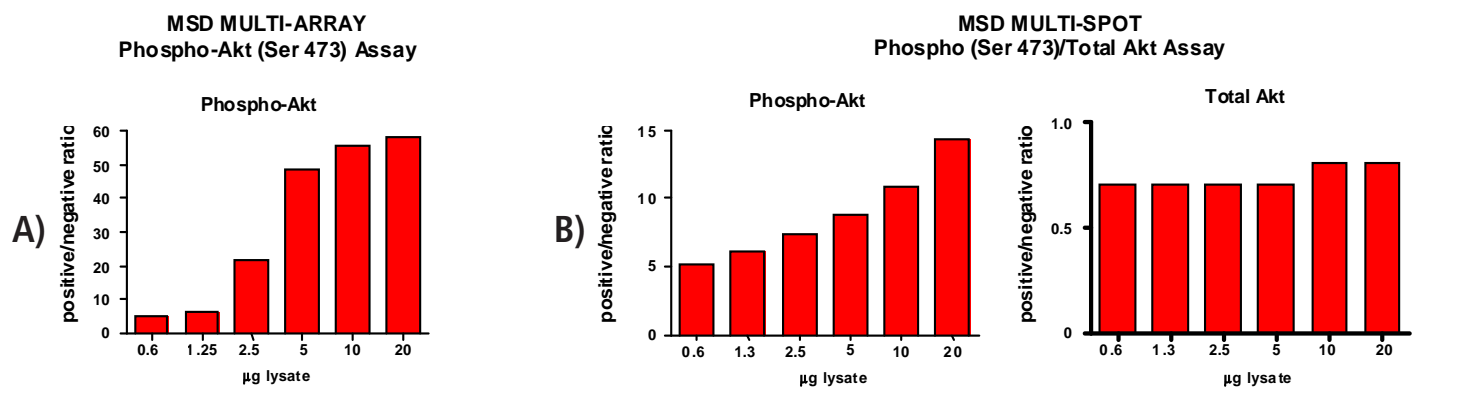
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-Akt (Ser 473)

Catalog No:	C11CA-1
Contents:	2 x 100 µg MSDLY0004 pAkt Negative Control Cell Lysate Cell lysate from growing Jurkat cells treated with 50 µM LY294002 for 2.5 hours to inhibit Akt phosphorylation 2 x 100 µg MSDLY0001 pAkt Positive Control Cell Lysate Cell lysate from growing Jurkat cells
Concentration:	2 mg/mL in MSD® Complete Tris Lysis Buffer
Volume:	2 vials (50 µL) negative lysate 2 vials (50 µL) positive lysate
Preparation:	Following cell treatment, Jurkat cell lysates were prepared on ice in MSD Complete Tris Lysis Buffer. Cell debris was cleared by centrifugation.
Storage:	Lysates should be stored at -80°C. Lysates will retain approximately 90% of activity after a single round of freeze thaw if handled properly (thawed on ice and immediately refrozen in smaller aliquots).
Quality Control:	Lysates have been tested for performance in Western Blot and MSD Assays.

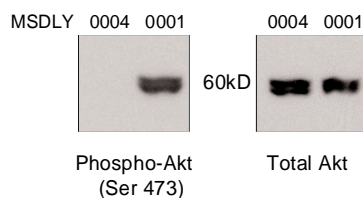
MSD MULTI-SPOT® Assay Results

Figures A and B (below) illustrate typical titrations for MSDLY0004 (pAkt negative) and MSDLY0001 (pAkt positive) cell lysates. Figure A depicts the MSD MULTI-ARRAY® Phospho-Akt (Ser 473) Assay; Figure B depicts the MSD MULTI-SPOT Phospho (Ser 473)/ Total Akt Assay. The results are presented as a ratio of the signals obtained with pAkt positive and pAkt negative lysates. In both assays, the phospho-Akt signal ratios increase with the amount of lysates. The signal ratios for total Akt in the Phospho/Total Akt Assay remain at a constant level close to one throughout the titration. The representative results shown below are for demonstration purposes only and individual results may vary depending upon experimental application.



Traditional Western Blot Results

MSDLY0004 and MSDLY0001 whole cell lysates (20 µg each) were analyzed by Western Blot with Phospho (Ser 473)-specific and Total Akt antibodies.



FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC OR THERAPEUTIC PROCEDURES.

20192-v1-2010Feb